

# 15-Ketoeicosa- 6E,8Z,11Z,13E- tetraenoic acid

Biosynthesis *via* 15-OH prostaglandin dehydrogenase from lung and liver microsomes and by a nonenzymatic process.

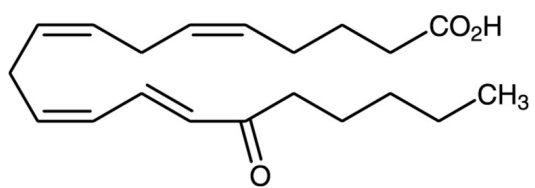
## Ordering Information

[Order Online »](#)

BML-K015-0050	50µg
---------------	------

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -80°C for up to 1 year. Product is subject to oxidation. Protect from air. Open ampule and aliquot into suitable tubes or vials. For experiments with cells or tissues, aliquots of ethanol solutions may be added directly to aqueous media. For HPLC, aliquots of ethanol solutions should be evaporated to dryness under a stream of nitrogen and HPLC mobile phase added to the desired concentration.
<b>Long Term Storage</b>	-80°C
<b>Shipping</b>	Dry Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Appearance</b>	Liquid.
<b>CAS</b>	81416-72-0
<b>MW</b>	318.5
<b>Purity</b>	≥95% (HPLC)



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)